

NPPB

Cl⁻ channel blocker

A potent chloride channel blocker (IC_{50} =0.1-100 μ M depending on channel subtype and assay method). Inhibits cyclooxygenase (COX; IC_{50} =8 μ M). Evidence of mitochondrial uncoupling in mouse peritoneal macrophages, due to protonophore activity.

Citations: 7

[View Online »](#)

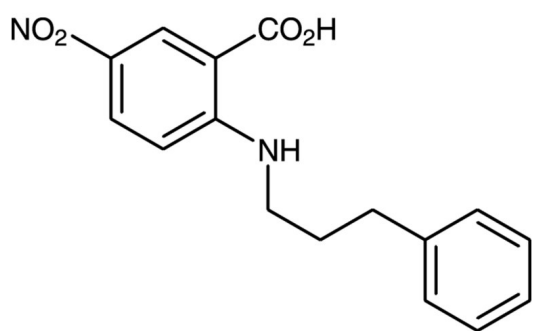
Ordering Information

[Order Online »](#)

BML-CM100-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at room temperature. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	5-Nitro-2-(3'-phenylpropyl-amino)benzoic acid
Appearance	Yellow solid.
CAS	107254-86-4
Couple Target	Chloride channel, COX
Couple Type	Blocker, Inhibitor
Formula	$C_{16}H_{16}N_2O_4$
MW	300.3
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (25mg/ml) or 100% ethanol (10mg/ml; warm).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com